

Macrophage marker monoclonal antibody (HAM56) (Ready-to-Use)

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 7

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-30935	~6 µg - 6ml
-----------	-------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application IHC (PS)

Clone HAM56

Crossreactivity Cross-reacts with a subpopulation of endothelial cells including those of capillaries and smaller blood vessels.

Formulation Liquid. *Ready-to-use* solution in PBS containing 0.1% sodium azide and 1% BSA.

Host Mouse

Immunogen Sonicated human alveolar macrophages.

Isotype IgM

Purity Detail Affinity purified.

Recommendation
Dilutions/Conditions Immunohistochemistry (paraffin sections; this antibody can be used for staining of paraffin-embedded tissue sections fixed in neutral buffered formalin and is compatible with most commonly used histological fixatives. Choose a good quality enzyme-linked anti-mouse IgM antibody for detection). Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually.

Source Purified from hybridoma tissue culture supernatant.

Species Reactivity Human, Monkey

Specificity Recognizes fixed macrophages such as “tingible macrophages”, interdigitating macrophages of lymph nodes or tissue macrophages.

Worry-free Guarantee



ENZO LIFE SCIENCES
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

ENZO LIFE SCIENCES
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

ENZO LIFE SCIENCES
France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

ENZO LIFE SCIENCES
Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com

This antibody is covered by our Worry-Free Guarantee